
DCAMKL1

Product code: 02-139

Full-length human DCAMKL1 [1-729(end) amino acids of accession number NP_004725.1] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 230 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 200
IC50 at 1 mM ATP (nM)	: 410
